

Separation of digested proteins on Silica gel 60

Plate	ProteoChrom® HPTLC Silica gel 60 F254s (P/N 1. 05650)
Mobile phase (v/v)	2-butanol/pyridin/ammonia25%/water (39/34/10/26)
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Detection	White light
Detection	0,5% Ninhydrin spray 20s and heat 1-2min at 110°C

Chromatographic Data

No.	compounds	Concentration (mg/ml)	Solvent	Application volume μ l
1	lysozyme digested	2	water/ACN 85/15	7
2	myoglobin digested	2	water/ACN 85/15	7
3	cytochrome c digested	2	water/ACN 85/15	7
4	β -casein digested	2	water/ACN 85/15	7
5	BSA digested	2	water/ACN 85/15	7

